

EED FL (GST)

CATALOG NO.: RD-11-498

LOT NO.:

DESCRIPTION: Human recombinant EED (2-441; NM_003797; 77.4 kDa) expressed in *E. coli* with an N-terminal GST-tag. EED functions as H3K27me3 reader within the PRC2/EED-EZH2 complex and mediates protein-protein interactions. During development, Polycomb Repressive Complex 2 (PRC2) is the principal methyltransferase responsible for generating trimethylated histone H3 lysine-27 (H3K27me3), an epigenetic mark essential for programmed repression of gene expression¹⁻⁵.

PURITY: >90% by SDS-PAGE

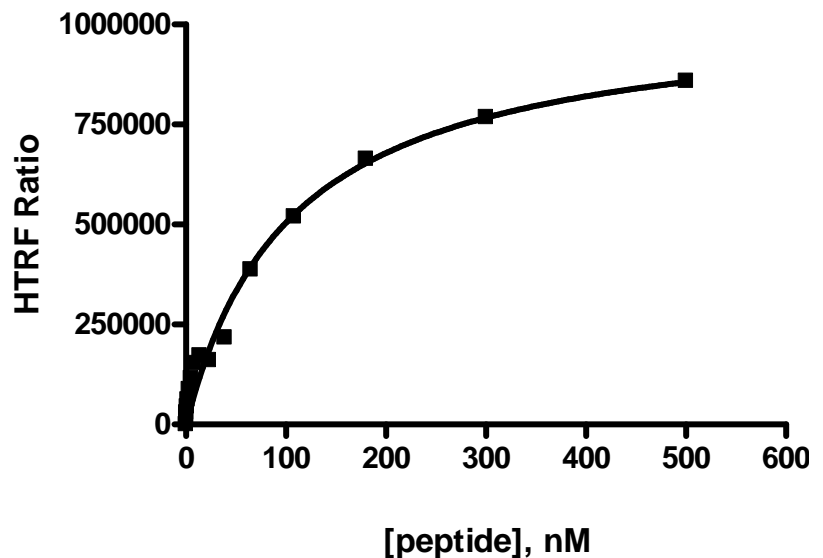
SUPPLIED AS: ____ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v)

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted protein is not recommended.

REFERENCES: 1) R. Cao *et al. Science* 2002 **298** 1039; 2) K. Plath *et al. Science* 2003 **300** 131; 3) J. Silva *et al. Dev. Cell* 2003 **4** 481; 4) S. Erhardt *et al. Development* 2003 **130** 4235; 5) R. Cao & Y. Zhang *Curr. Opin. Genet. Dev.* 2004 **14** 155



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4 µg of RBC EED FL (GST). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



EED/peptide interaction Assay. EED FL binding to B-H3(21-44) peptide was assessed using Cisbio HTRF detection (cat #61GSTTLB and 610SAXLB). The 25µL reaction contained 5nM protein, variable concentration of peptide and detection mix. Fluorescence emission (665 and 615nm) was read using Envision reader (Perkin Elmer) following 1.5h incubation.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.